

# PRODUCT INFORMATION



## STING M284 variant (human, recombinant)

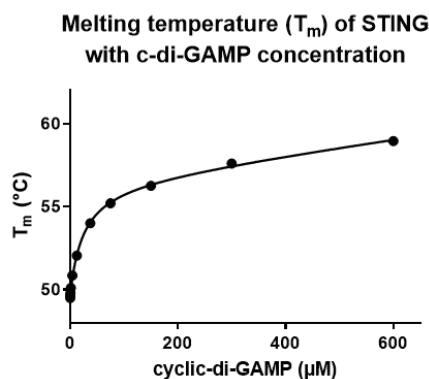
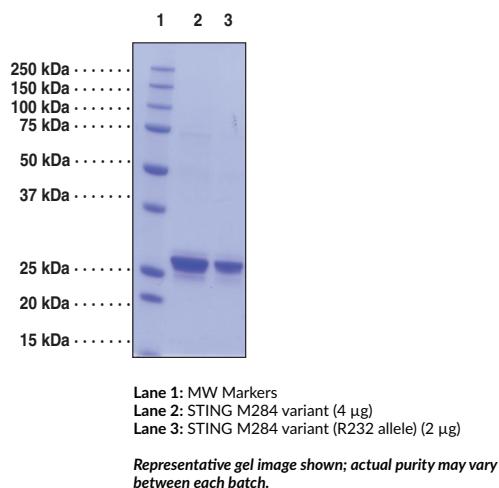
Item No. 23594

### Overview and Properties

<b>Synonyms:</b>	Endoplasmic Reticulum Interferon Stimulator, ERIS, hSTING, MITA, Stimulator of Interferon Genes Protein, TMEM173, Transmembrane Protein 173
<b>Source:</b>	N-terminal Histidine-tagged human recombinant protein expressed in <i>E. coli</i>
<b>Amino Acids:</b>	138-379 (N-terminal truncation)
<b>Uniprot No.:</b>	Q86WV6
<b>Molecular Weight:</b>	28.8 kDa
<b>Storage:</b>	-80°C (as supplied); avoid freeze/thaw cycles by storing protein in aliquots
<b>Stability:</b>	≥2 years
<b>Purity:</b>	<i>batch specific</i> (≥80% estimated by SDS-PAGE)
<b>Supplied in:</b>	50 mM HEPES, pH 8.0 with 150 mM sodium chloride and 10% glycerol
<b>Protein Concentration:</b>	<i>batch specific</i> mg/ml
<b>Activity:</b>	Serial dilutions of canonical 3'3'-cGAMP were incubated with 5 µg recombinant human STING M284 variant in 50 mM HEPES, pH 7.5, 150 mM sodium chloride, 10% glycerol, and SYPRO® Orange dye at 4°C. <sup>1</sup> The reaction was read on a BioRad CFX96 Touch™ Real-Time PCR Detection System at 4-100°C. <sup>2</sup> The binding of the ligand stabilizes the protein structure, increasing the melting temperature ( $T_m$ ), which is detected via a thermal shift assay (TSA), also known as a differential scanning fluorimetry (DSF) assay. <sup>2</sup>

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Images



Binding Activity of STING M284 variant (Item No. 24594). STING M284 variant (5 µg) was incubated with serial dilutions of 3'3'-cGAMP (Item No. 17966) and SYPRO® Orange dye. The detected increase in  $T_m$  indicates binding.

WARNING  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/24/2020

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

STING M284 variant (human, recombinant) contains amino acids 138-379 of the wild-type variant (R232) with a methionine substituted for arginine at position 284. Stimulator of interferon genes (STING) is a component of the innate immune response that binds to cyclic dinucleotides, which are bacterial second messengers, leading to activation of NF- $\kappa$ B and transcription of immunomodulatory genes, including type I interferon (IFN).<sup>3-6</sup> The R232 variant of STING is the most common variant in the human population, found at a frequency of 57.9% in the 1000 Genome Project.<sup>7</sup> The SNP variant H232 is found at a 13.7% frequency. The R284M mutation in STING is associated with constitutive activation of downstream signaling. It increases the propensity of STING to dimerize and associate with the kinase TBK1, enhancing the ability of STING to activate IRF3 and NF- $\kappa$ B and induce a type I IFN response.<sup>8</sup> However, the R284M mutation occurs outside of the dimerization region between positions 153-177, so rather than a direct effect on dimerization, it is predicted to promote or inhibit binding of a cellular factor that stabilizes or impairs STING dimerization.

## References

---

1. Diner, E.J., Burdette, D.L., Wilson, S.C., *et al.* The innate immune DNA sensor cGAS produces a noncanonical cyclic dinucleotide that activates human STING. *Cell Rep.* **3(5)**, 1355-1361 (2013).
2. Niesen, F.H., Berglund, H., and Vedadi, M. The use of differential scanning fluorimetry to detect ligand interactions that promote protein stability. *Nat. Protoc.* **2(9)**, 2212-2221 (2007).
3. Sun, L., Wu, J., Du, F., *et al.* Cyclic GMP-AMP synthase is a cytosolic DNA sensor that activates the type I interferon pathway. *Science* **339(6121)**, 786-791 (2013).
4. Wu, J., Sun, L., Chen, X., *et al.* Cyclic GMP-AMP is an endogenous second messenger in innate immune signaling by cytosolic DNA. *Science* **339(6121)**, 826-830 (2013).
5. Konno, H., Konno, K., and Barber, G.N. Cyclic dinucleotides trigger ULK1 (ATG1) phosphorylation of STING to prevent sustained innate immune signaling. *Cell* **155(3)**, 688-698 (2013).
6. Burdette, D.L., Monroe, K.M., Sotelo-Troha, K., *et al.* STING is a direct innate immune sensor of cyclic-di-GMP. *Nature* **478(7370)**, 515-518 (2011).
7. Yi, G., Brendel, V.P., Shu, C., *et al.* Single nucleotide polymorphisms of human STING can affect innate immune response to cyclic dinucleotides. *PLoS One* **8(10)**, e77846 (2013).
8. Tang, E.D. and Wang, C.-Y. Single amino acid change in STING leads to constitutive active signaling. *PLoS One* **10(3)**, e0120090 (2015).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM